

p53 (human), (recombinant) (GST- tag)

p53 is a much studied and complex multifunctional protein, which acts as a tumor suppressor in many tumor types and induces growth arrest or apoptosis depending on the physiological circumstances and cell type. It is involved in cell cycle regulation as a trans-activator that acts to negatively regulate cell division by controlling a set of genes required for this process. One of the activated genes is an inhibitor of cyclin-dependent kinases. Apoptosis induction appears to be mediated either by stimulation of BAX and FAS antigen expression, or by repression of Bcl-2 expression. The p53 pathway responds to stresses that can disrupt the fidelity of DNA replication and cell division. A stress signal is transmitted to the p53 protein by post-translational modifications. Such modification results in the activation of p53 protein as a transcription factor that initiates a program of cell cycle arrest, cellular senescence or apoptosis.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

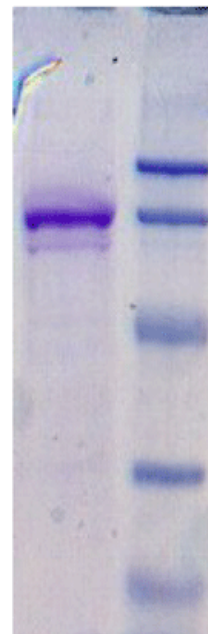
BML-FW9370-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot of prod. no. BML-FW9370 after SDS-PAGE followed by blotting onto PVDF and probing with mouse monoclonal antibody to p53 (Cat. #BML-SA293, 1:5000).



1 2

Coomassie stained gel of prod.no. BML-FW9370: Lane 1: FW9370 (3 μ g), Lane 2: MW markers (top to bottom) 113, 93, 50, 35, and 28 kDa.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TRP53, Tumor protein p53

Formulation Liquid. In 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM EGTA, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol.

MW ~70kDa (~85/90kDa observed with GST tag)

Purity Detail Purified using glutathione sepharose chromatography.

Source Produced in *E. coli*. Full length human p53 is fused to a GST-tag.

UniProt ID P04637



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)